

Reference

1. Engvall, E. and Perlmann, P., *Immunochemistry*, **8**, 871 (1971).



Avidin-Alkaline Phosphatase

**Catalog Number
170-6533**



Specifications

Contents	1 ml
Preparation of conjugate	Avidin is conjugated to alkaline phosphatase using a modification of the one step glutaraldehyde procedure described by Engvall and Perlmann. ¹
Buffer	10 mM Tris, 150 mM sodium chloride, 1.0 mM magnesium chloride, pH 8.0. Contains 1.0% bovine serum albumin as a stabilizer and 0.1% sodium azide as a bacteriostat.
Storage conditions	Avidin-AP will arrive frozen. It may be stored frozen until needed. Once thawed, store at 4 °C. Repeated freeze-thaw cycles should be avoided.
Shelf life	1 year at 4 °C.

Recommended working dilution 1:1,000 when used with NHS-biotinylated proteins. 1:3,000 when used with the biotinylated standards. Different dilutions may be optimal for other applications.

Additional Information

The instruction manuals provided with the Biotinylated SDS-PAGE Molecular Weight Standards Avidin-AP Kits (catalog numbers 161-0308, 161-0313, and 161-0322) contain complete details for assay procedures. To biotinylate proteins use Bio-Rad's NHS-Biotin (catalog number 170-6529).